

# CD8B, Fc Fusion, Avi-Tag, Biotin-Labeled Recombinant

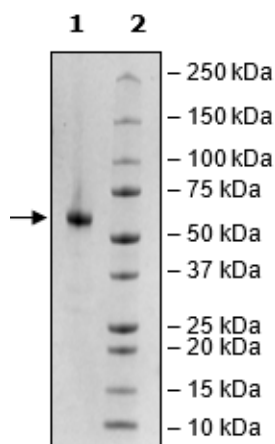
Catalog: 101842  
Lot: 230725

## Product Information

<b>Description:</b>	Recombinant human CD8b (cluster of differentiation 8 b), encompassing amino acids 22-170. This construct contains a C-terminus Fc region of human IgG1 followed by an Avi-Tag™. This protein was affinity purified.
<b>Background:</b>	CD8 is a co-receptor for the TCR (T cell receptor) in T cells, binding to MHC (major histocompatibility complex) class I proteins. CD8 is a typical marker of cytotoxic T cells and is involved in signaling. CD8 has two isoforms, a and b. CD8b recruits Lck (lymphocyte-specific protein tyrosine kinase) to the TCR-CD3 complex, and Lck phosphorylates multiple proteins involved in activation of cytotoxic T-lymphocytes. It is thus critical for the lysis of cancer cells.
<b>Species:</b>	Human
<b>Construct:</b>	CD8B (22-170-Fc(IgG1)-Avi)-(Biotin)
<b>Concentration:</b>	1.02 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	46 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_172213.5
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful for SDS-PAGE and avidin pulldown assays.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

